

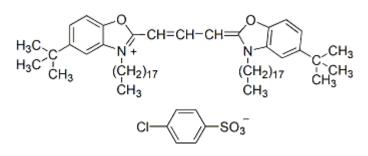
Neuro-DiO

Biotium developed Neuro-DiO as a replacement for the green fluorescent DiO, which has been found to be difficult to use for neurons and cell suspension due to low solubility, tendency to form aggregates and slow lateral diffusion rate.

Product Description

Biotium developed Neuro-DiO as a replacement for the green fluorescent DiO, which has been found to be difficult to use for neurons and cell suspension due to low solubility, tendency to form aggregates and slow lateral diffusion rate. Neuro-DiO and DiO have nearly identical absorption and emission spectra, but the former has better solubility in membranes and does not form nonfluorescent aggregates, which also tend to slow down the dye diffusion rate in membranes. We also offer Neuro-DiO in Vegetable Oil for microinjection studies, and in our CellBrite™ Green Cytoplasmic Membrane Dye, a ready-to-use dye solution for cell labeling.

- $\lambda_{Ex}/\lambda_{Em} = 484/501 \text{ nm}$
- Yellow-orange solid soluble at 1-2 mM (with heating) in DMF, DMSO, or ethanol
- Soluble at 1-2 mM in vegetable oil with heating and sonication
- Store at 4°C, protect from light
- C67H105CIN2O5S
- MW: 1086



This datasheet was generated on November 23, 2025 at 03:08:31 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/neuro-dio/

Product attributes

Call us: 800-304-5357

Probe cellular localization	Membrane/cell surface, Membrane/vesicular
For live or fixed cells	For fixed cells, For live/intact cells
Assay type/options	Co-cultures, Extended staining (several days to weeks)
Fixation options	Fix before staining (formaldehyde), Fix after staining (formaldehyde), Permeabilize before staining
Colors	Green
Excitation/Emission	484/501 nm